

Tecan Schedules Analysts' and Media Conference on March 4 to Discuss Full-Year 2009 Financial Results

Männedorf, Switzerland, February 26, 2010 – The Tecan Group AG (SIX Swiss Exchange: TECN) will hold an Analysts' and Media Conference to discuss the full-year 2009 financial results on Thursday, March 4, 2010 at 10:00 a.m. CET. The results will be published on March 4, 2010 at 7:00 a.m. CET.

The conference will be audio-webcast with synchronized presentation slides. Interested parties may access the webcast via Tecan's homepage www.tecan.com. A link to the webcast will be provided immediately prior to the event.

The conference can also be heard by phone.

The dial-in numbers for the call are as follows:

Participants from Europe: +41 91 610 5600 or +44 207 107 0611 (UK)

Participants from the U.S.: +1 866 291 4166

Please dial in 15 minutes before the beginning of the meeting.

The recorded webcast will be available the same day on the website www.tecan.com together with the media release, the presentation and a preprint of the Financial Report 2009.

About Tecan

Tecan (www.tecan.com) is a leading global provider of laboratory instruments and solutions in biopharmaceuticals, forensics, and clinical diagnostics. The company specializes in the development, production and distribution of automation solutions for laboratories in the life sciences sector. With its subsidiary REMP (www.remp.com), Tecan is the market leader in automated laboratory storage and logistics systems. Its clients include pharmaceutical and biotechnology companies, university research departments, forensic and diagnostic laboratories. As an original equipment manufacturer, Tecan is also a leader in developing and manufacturing OEM instruments and components that are then distributed by partner companies. Founded in Switzerland in 1980, the company has manufacturing, research and development sites in both Europe and North America and maintains a sales and service network in 52 countries.

In 2008, Tecan generated sales of CHF 396 million (USD 366.7 million; EUR 250.7 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TK: TECN/Reuters: TECZn.S/ ISIN CH0012100191).



2

For further information, please contact:

Tecan Group
Martin Braendle
Head of Corporate Communications &
Investor Relations
Tel. +41 (0) 44 922 84 30
Fax +41 (0) 44 922 88 89
investor@tecan.com
www.tecan.com